

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IFN- γ Kit
Kit Catalog Numbers: K152QO-Series
Kit Lot Number: K0081505
Expiration Date: 31 Aug 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 140%	< 25%	90% - 140%	< 25%	90% - 140%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IFN- γ	Calibrator	764	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	469	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	32	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	2.12	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
764	521,000	1.4
191	131,000	1.5
47.8	32,400	1.6
11.9	8,270	1.7
2.98	2,170	1.4
0.746	594	1.8
0.187	192	2.9
0.00	66	4.4

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0047351	2-8°C	31 Dec 2022
Anti-ms IFN- γ Antibody	D0081378	2-8°C	31 Mar 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080250	2-8°C	28 Feb 2025
Diluent 41	M0410033	\leq -10°C	31 Oct 2022
Diluent 45	M0450033	\leq -10°C	31 Aug 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 AUG 2020